
MAP3K1

Product code: 07-103

Human MAP3K1, catalytic domain [1193-1512(end) amino acids of accession number NP_005912.1] was expressed as N-terminal GST-fusion protein (62 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : MAP2K1
Cascade Assay *
ATP (μ M) : 1000
Metal : Mg
Reference compound : Staurosporine
IC50 at 1 mM ATP (nM) : 160

* MAP2K1/Erk2/Modified Erktide
